

PRODUCT DATA SHEET

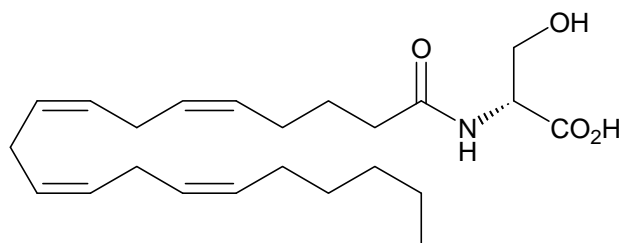


PRODUCT: N-Arachidonoyl-L-serine

CATALOG NO.: FA-037

LOT NO.:

STRUCTURE:



PHYSICAL APPEARANCE: Waxy solid

MOLECULAR FORMULA: C₂₃H₃₇NO₄

MOLECULAR WEIGHT: 391.5

PURITY: 98% (TLC: 30% Methanol/methylene chloride; R_f = 0.17)

SOLUBILITY: Soluble in DMSO (4 mg/mL)

STORAGE: Store, as supplied, at -80°C for up to 1 year. Store solutions at -80°C for up to 1 month.

APPLICATION NOTES: N-Arachidonoyl-L-serine is a recently discovered endocannabinoid isolated from bovine brain. It binds only very weakly to cannabinoid CB1 and CB2 receptors as well as vanilloid TRPV1 receptors. However, it produces endothelium-dependent vasodilation in rat isolated mesenteric arteries and abdominal aortas. It also suppresses LPS-induced TNF α production in a murine macrophage cell line and in mice. Its biological profile closely parallels that of abnormal cannabidiol, an agonist at a putative novel cannabinoid-like receptor. Thus, N-arachidonoyl-L-serine may be an endogenous agonist for this receptor.

REFERENCES: 1. G. Milman *et al. Proc. Natl. Acad. Sci. USA* 2006 **103** 2428

The pharmacological and toxicological properties of this product have not been fully investigated. Exercise caution in use and handling. This product must not be used in humans.